



B-Phycoerythrin Certificate of Analysis

Product Number D4-000
Concentration 24.18 mg/ml
Store at 2-8°C

Form/Shipping & Storage

Supplied as a functionally water soluble ammonium sulfate precipitate. Upon receipt, store at 2-8°C in the dark, do not freeze. Phycobiliproteins are sensitive to freeze-thaw cycles.

Handling

Avoid exposure to heat and light.

Buffer

60% ammonium sulfate in 100 mM phosphate and 100 mM Sodium Chloride buffer pH 7.4 with 2mM Sodium Azide.

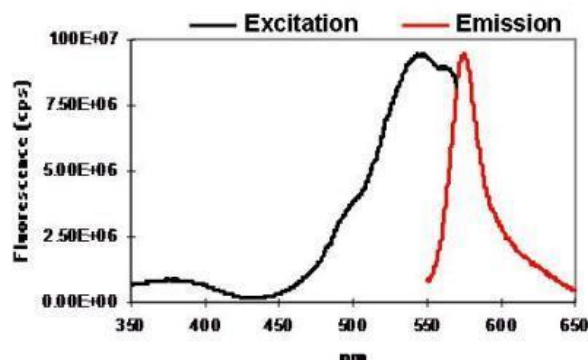
Note

For research use only, not for diagnostic or therapeutic use.

Method

The material concentration was measured using a dilution of well mixed material in 100 mM phosphate buffer pH 7.4. The reported result is the average over all dilutions and all readings.

The HPLC was run using 25 µl injections in 10 mM phosphate buffer pH 7.4 with 150 mM sodium chloride and 2 mM sodium azide.



The fluorometer run was performed by diluting 300 µl of 0.01 mg/mL B-PE in 3 mL of 100 mM. phosphate buffer pH 7.4 with 100 mM sodium chloride and 2 mM sodium azide.

Specifications:	Actual Result:
$A_{545} / A_{280} :$	≥ 5.50 5.92
$A_{620} / A_{545} :$	≤ 0.01 0.0019
$A_{556} / A_{495} :$	≤ 2.50 2.43
> 95% single peak by HPLC	100.0%
Ex at 532 nm peak 565-585	573.0 nm

4985 Winchester Blvd. Frederick, MD 21703
301.732.5415

www.columbiabiosciences.com

For technical inquiries: info@columbiabiosciences.com
For sales inquiries: sales@columbiabiosciences.com